

Complement component C1s (human) monoclonal antibody (9)

C1s proenzyme is a 79.8kDa glycoprotein produced primarily by hepatocytes but also by other cells. The proenzyme is produced as a single-chain polypeptide of 688 amino acids. Upon activation, the peptide is cleaved between R437 and I438, producing an activated enzyme consisting of two disulfide-linked chains, A chain (52.2kDa) and B chain (27.7kDa). The calcium-dependent catalytic subunit of the C1 complex (C1q, C1r, C1s) is comprised of two C1s and two C1r molecules, and this serine protease tetramer is responsible for the initial steps in the cleavage of C4 and C2 in the classical complement pathway.

This antibody is covered by our [Worry-Free Guarantee](#).

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Long Term Storage +4°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name	C1 esterase, Complement C1s subcomponent, C1s proenzyme
Application	ELISA, IHC
Clone	9
Formulation	Liquid. In 0.01M phosphate buffer, pH 7.4, containing 0.5M sodium chloride and 15mM sodium azide.
GenBank ID	716
Host	Mouse
Immunogen	Human plasma complement component C1s Proenzyme.
Isotype	IgG1κ
Purity Detail	Protein A/G-purified.
Source	From culture supernatant.
Species Reactivity	Human
Specificity	Specific for C1s proenzyme and activated free C1s.
UniProt ID	P09871

Worry-free Guarantee

This antibody is covered by our [Worry-Free Guarantee](#)



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
[info-
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
[info-
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France
Phone: +33 472 440 655
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
[info-
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)

Last modified: May 29, 2024